



6405 Mira Mesa Blvd, Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Data Sheet

ANPRA/Aequorin Expression Vector Catalog #: 60641

Description: The ANPRA/Aequorin expression vector is designed to co-express human atrial natriuretic peptide receptor A (ANPRA, also called natriuretic peptide receptor A/guanylate cyclase A) and jellyfish (*Aequorea victoria*) Aequorin in mammalian cells.

Application: Suitable for use in transient assays with the PDE1B-HEK293 Cell Line (BPS Bioscience #60402).

Size: 20 µg

Concentration: 80 ng/µl

Storage buffer: 10 mM Tris-HCl (pH 8), 1 mM EDTA

Storage condition: Store plasmid at -20°C in aliquots. Avoid repeated freeze/thaw cycles.

Stability: Stable for at least one year when stored as recommended.

Note: The vector is suitable for transient transfection. It is not suitable for amplification in bacteria.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com